

PRODUCT INFORMATION



Hyperforin (dicyclohexylammonium salt)

Item No. 19572

CAS Registry No.: 238074-03-8

Formal Name: (1R,5S,6R,7S)-4-hydroxy-6-methyl-1,3,7-tris(3-methyl-2-buten-1-yl)-5-(2-methyl-1-oxopropyl)-6-(4-methyl-3-penten-1-yl)-bicyclo[3.3.1]non-3-ene-2,9-dione, N-cyclohexylcyclohexanamine (1:1)

Synonyms: Hyp-DCHA, Hyperforin-DCHA

MF: $C_{35}H_{52}O_4 \cdot C_{12}H_{23}N$

FW: 718.1

Purity: $\geq 95\%$

UV/Vis.: λ_{max} : 297 nm

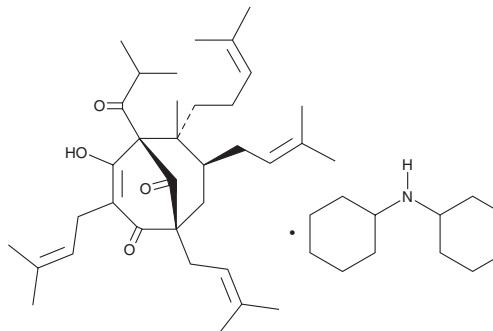
Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years

Special Conditions: Keep under inert gas. Protect from light.

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.



Laboratory Procedures

Hyperforin (dicyclohexylammonium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the hyperforin (dicyclohexylammonium salt) in the solvent of choice, which should be purged with an inert gas. Hyperforin (dicyclohexylammonium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of hyperforin (dicyclohexylammonium salt) in these solvents is approximately 20 mg/ml.

Hyperforin (dicyclohexylammonium salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, hyperforin (dicyclohexylammonium salt) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Hyperforin (dicyclohexylammonium salt) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Hyperforin is a natural activator of the steroid X receptor (active at 0.1 to 0.5 $\mu g/ml$) and inhibitor of several cytochrome P450 (CYP) isoforms ($IC_{50} = 10 \mu g/ml$ for CYP2D6).^{1,2} It can also inhibit microsomal prostaglandin E_2 synthase-1 (mPGES-1; $IC_{50} = 1 \mu M$), 5-lipoxygenase activating protein, and sirtuins.³⁻⁵

References

1. Wentworth, J.M., Agostini, M., Love, J., et al. *J. Endocrinol.* **166**, R11-R16 (2000).
2. Obach, R.S. *J. Pharmacol. Exp. Ther.* **294**(1), 88-95 (2000).
3. Chang, H.H. and Meuillet, E.J. *Future Med. Chem.* **3**(15), 1909-1934 (2011).
4. Greiner, C., Hörnig, C., Rossi, A., et al. *Br. J. Pharmacol.* (2011).
5. Milne, J.C. and Denu, J.M. *Curr. Opin. Chem. Biol.* **12**, 11-17 (2008).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/26/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM